

VivaFix™ Cell Viability Assays



Bio-Rad's VivaFix Cell Viability Assays provide a sensitive measurement for determining the viability of mammalian cells by flow cytometry and microscopy.

Key Benefits of VivaFix Cell Viability Assays

- Optimal discrimination between dead and live cell populations
- Compatible with cell fixation
- 8 different excitation/emission wavelength combinations to fit the most demanding multicolor experiments

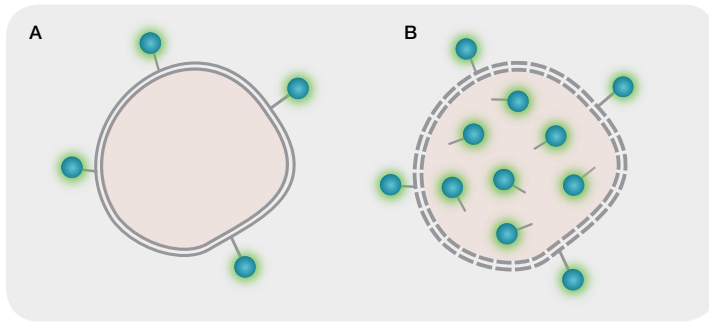
Visit bio-rad.com/web/vivafix for more information.

BIO-RAD

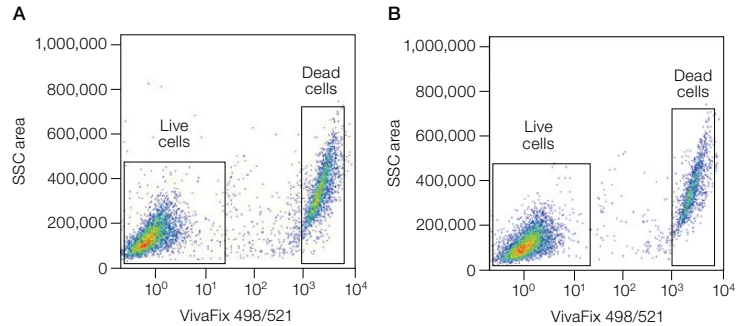
Accurately Determine the Viability of Your Cells

Taking advantage of an array of proprietary amine reactive dyes, VivaFix Cell Viability Assays can easily assist researchers with distinguishing between live and dead cells, providing at least a 100-fold difference in fluorescence intensity between the two populations. In addition, the brightness of VivaFix dyes is preserved upon treatment with fixative agents and can be used with biohazardous samples to track cell viability after fixation.

VivaFix Assays are available in a wide range of excitation and emission spectra and can be easily combined into any multicolor experiment. Choose between eight different assays to select the most appropriate excitation/emission wavelength for your experiments.



VivaFix Cell Viability Assay chemistry. **A**, VivaFix dyes bind to the cell surface primary amines of viable cells; **B**, in dead cells, where the plasma membrane is compromised, VivaFix dyes are able to permeate the cell and also react with intracellular primary amines. As a result, a greater number of fluorophores is associated with dead cells and at least a 100-fold difference in fluorescence intensity is measured between the live and the dead cells, thereby allowing an easier discrimination between the two populations.



Excellent separation between live and dead cells using the VivaFix Cell Viability Assay. Jurkat cells were stained with the VivaFix 498/521 dye, fixed with 3.7% formaldehyde (**A**) or not fixed (**B**), and analyzed with the S3™ Cell Sorter. SSC, side scatter.

Ordering Information

Catalog Number	VivaFix Cell Viability Assay	Excitation Maximum, nm	Emission Maximum, nm	Optimal Excitation Laser, nm	Optimal Filter Set for S3 Cell Sorter
135-1111	VivaFix 353/442	353	442	355	—
135-1112	VivaFix 410/450	410	450	405	—
135-1113	VivaFix 408/512	408	512	405	—
135-1114	VivaFix 398/550	398	550	405	—
135-1115	VivaFix 498/521	498	521	488	FL1 (525/30)
135-1116	VivaFix 547/573	547	573	561	FL2 (586/25)
135-1117	VivaFix 583/603	583	603	561	FL3 (615/25)
135-1118	VivaFix 649/660	649	660	640	—



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 3065 7550
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026 674 55 05
Taiwan 886 2 2578 7189 **Thailand** 1800 88 22 88 **United Kingdom** 020 8328 2000